

Anti-p21CIP1 (p21/WAF1)

Mouse Monoclonal Antibody

Catalog # **C27-61M**

Lot # Z2014-21

Cited Applications

ELISA, WB, IP

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the p21CIP1 (p21/WAF1) protein

Cross Reactivity

Human

Host/Isotype/Clone#

Mouse, IgG

Immunogen

Produced by repeated immunizations with recombinant human p21 fusion protein.

Formulation

0.02 M Potassium Phosphate, 0.5 M Sodium Chloride, pH 7.2, 0.01% (w/v) Sodium Azide

Stability

1yr at -20°C from date of shipment

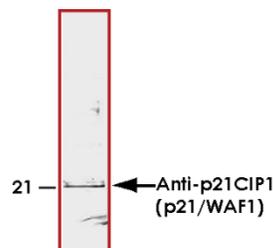
Scientific Background

CIP 1 (Cyclin-Dependent Kinase Inhibitor 1A) regulates cell cycle progression, terminal differentiation, and apoptosis (1). CIP1 is induced by p53 and is a potent inhibitor of cyclin-dependent kinase (CDK) activity. DNA damage leads to increased expression of CIP1 in cyclin E-containing complexes and to an associated decrease in cyclin-dependent kinase activity. CIP1 is a critical downstream effector in the p53-specific pathway of growth control in mammalian cells (2).

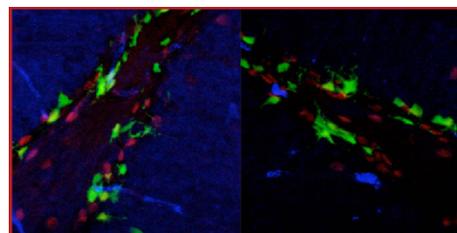
References

1. el-Deiry, W S. et al: WAF1/CIP1 is induced in p53-mediated G1 arrest and apoptosis. *Cancer Res.* 1994;54(5):1169-74.
2. Wagner, M. et al: Cyclin-dependent kinase-inhibitor 1 (CDKN1A) in the squamous epithelium of the oropharynx: possible implications of molecular biology and compartmentation. *Anticancer Res.* 2001; 21(1A):333-45.

Sample Data



Anti-p21CIP1 (p21/WAF1) detects human p21 by western blot. 10 µg of a HeLa whole cell lysate was loaded per lane and incubated with a 1:1,000 of Anti-p21CIP1 (p21/WAF1). A single band corresponding to human p21WAF1 is detected at ~21 kDa.



Immunohistochemistry of Anti-p21CIP1 (p21/WAF1) antibody 1:250. Tissue: human brain. Co-stained with YFP and Sox2 antibodies. Localization: p21 is nuclear and cytoplasmic. Staining: p21CIP1 as precipitated blue with Cy5, YFP as precipitated green with Cy2, and Sox2 as precipitated red with Cy3. z-stacks from confocal expressed as one composite focal plane.

Anti-p21CIP1 (p21/WAF1)

Mouse Monoclonal Antibody

Catalog #	C27-61M
Specific Lot #	Z2014-21
Purification	Affinity chromatography
Stability	1yr at -20°C from date of shipment
Storage & Shipping	Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.